

HAP1 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat HAP1

Instruction Manual

Catalog Number	PK-AB718-4205
Synonyms	HAP1 Antibody; Huntingtin-associated protein 1, neuroan 1, HLP, HIP5
Description	Huntington's disease (HD), a neurodegenerative disorder characterized by loss of striatal neurons, is caused by an expansion of a polyglutamine tract in the HD protein huntingtin. HAP1 was initially identified through a two-hybrid library screening; the binding of HAP1 to huntingtin correlated with the expansion of the polyglutamine tract. HAP1 also interacts with two cytoskeletal proteins (dynactin and pericentriolar autoantigen protein 1), suggesting that HAP1 may play a role in vesicular trafficking or organelle transport. HAP1 is also involved with the huntingtin-enhanced BDNF transport along the cellular microtubules. Attenuation of this process led to the loss of neurotrophic support and neuronal toxicity, which suggests that loss of this function might contribute to pathogenesis. Several alternatively spliced isoforms have been described for HAP1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	HAP1 antibody was raised in rabbits against a 19 amino acid peptide from near the center of human HAP1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. HAP1 antibody can be used for detection of HAP1 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Borrell-Pagès M, Zala D, Humbert S, et al. Huntington's disease: from huntingtin function and dysfunction to therapeutic strategies. <i>Cell Mol. Life Sci.</i> 2006; 63:2642-60. Li X-J, L S-H, Sharp AH, et al. A huntingtin-associated protein enriched in brain with implications for pathology. <i>Nature</i> 1995; 378:398-402. Engelender S, Sharp AH, Colomer V, et al. Huntingtin-associated protein 1 (HAP1) interacts with the p150(Glued) subunit of dynactin. <i>Hum. Molec. Genet.</i> 1997; 6:2205-12. Gauthier LR, Charrin BC, Borrell-Pages M, et al. Huntingtin controls neurotrophic support and survival of neurons by enhancing BDNF vesicular transport along microtubules. <i>Cell</i> 2004; 118:127-38.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4205P; HAP1 Peptide Cat.No. PK-AB718-1403; Mouse Brain Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.